

TCPOBOP

Constitutive androstane receptor (CAR) activator

TCPOBOP is an activator of the mouse constitutive androstane receptor (CAR) and a potent inducer of cytochrome P450 monooxygenase activity. Has been shown to have a potent tumor-promoting activity and hepatocarcinogenic effects in mice but not in rats.

Citations: 4

[View Online »](#)

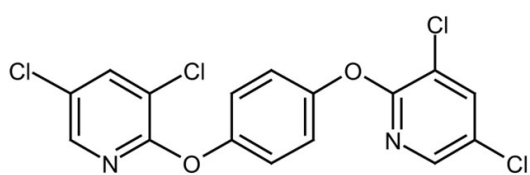
Ordering Information

[Order Online »](#)

BML-GR245-0025	25mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1,4-Bis-[2-(3,5-dichloropyridyloxy)]benzene

Appearance White solid.

CAS 76150-91-9

Couple Target Constitutive androstane receptor, Cytochrome P450

Couple Type Activator, Ligand

Formula $C_{16}H_8Cl_4N_2O_2$

MW 402.1

Purity $\geq 98\%$ (TLC)

Solubility Soluble in DMSO (5mg/ml).

Source Synthetic.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com